

Elham Hatami<sup>1\*</sup>, Thomas McCarthy<sup>2\*</sup>, Hye-Won Song<sup>1</sup>, Quyen Bao<sup>2</sup>, Youngsook Kim<sup>2</sup>, Hongduan Huang<sup>2</sup>, Rosary Nguyen<sup>2</sup>, Punya Narayan<sup>2</sup>, Zhiqi Zhang<sup>1</sup>, Samuel Shi<sup>1</sup>, Ruifang Li<sup>2</sup>, Larry Wang<sup>2</sup> and Aruna Ayer<sup>2</sup>

\*Authors contributed equally

<sup>1</sup>BD Biosciences, 10975 Torreyana Rd, San Diego, CA 92121, <sup>2</sup>BD Biosciences, 2350 Qume Drive, San Jose, CA, 95131

## Abstract

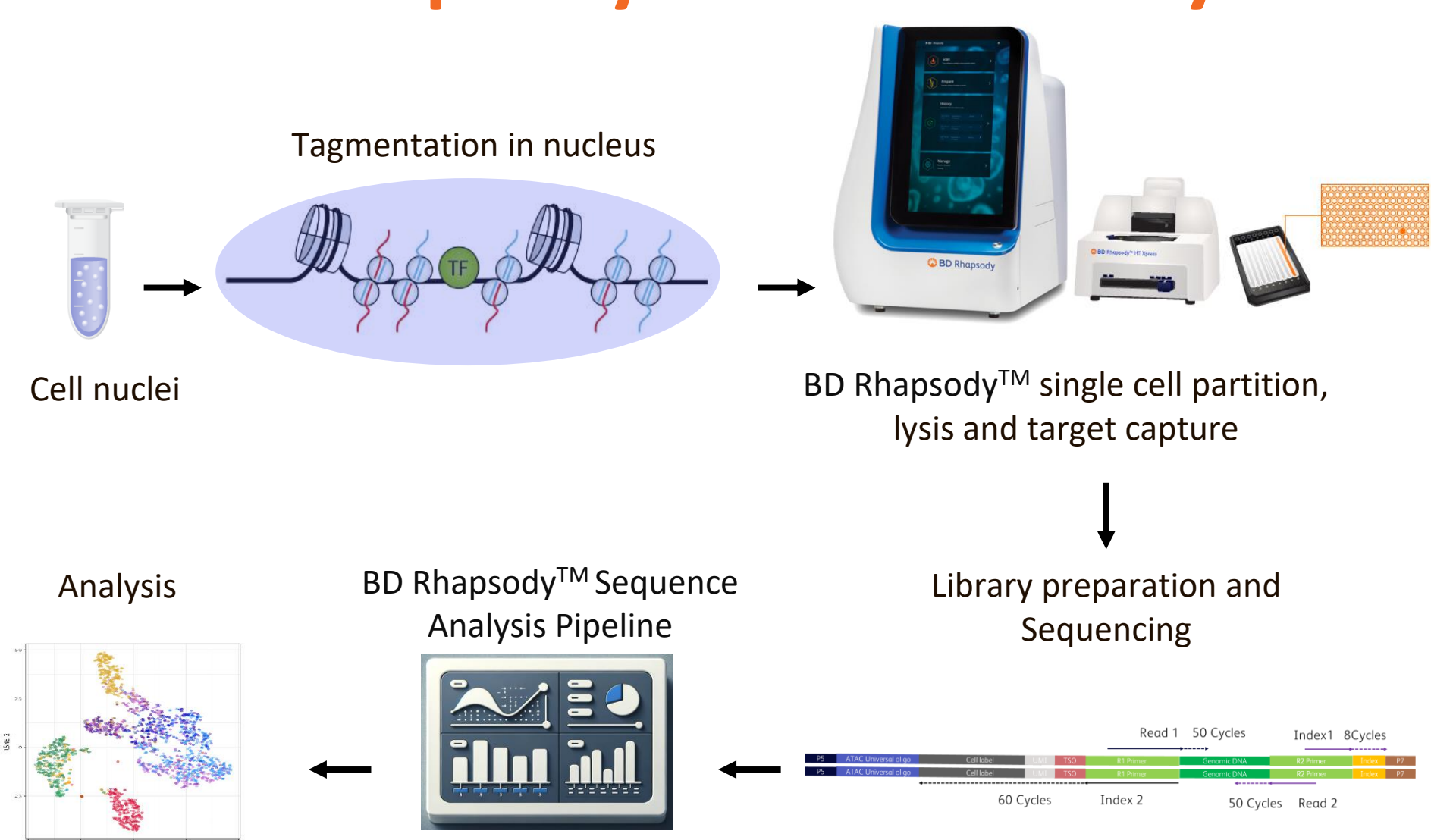
Single-cell RNA sequencing (scRNA-Seq) has significantly advanced our understanding of cell development and heterogeneity in complex systems at cellular resolution. However, its capacity to elucidate cell states and uncover gene regulatory programs remains limited. In contrast, chromatin state profiles serve as a valuable resource for gauging gene expression potential and provide insights regarding transcriptional regulation. When integrated with gene expression data, chromatin accessibility region (CAR) profiles offer a means to define the fundamental gene regulatory logic as the foundation of cell fate.

ATAC-Seq (Assay for Transposase-Accessible Chromatin using Sequencing) is a potent approach for profiling genome-wide CARs. To capture both mRNA gene expression data and gene regulatory information in a single assay, we conducted multiomic scATAC-Seq + scRNA-Seq on PBMCs, utilizing the gentle and robust microwell-based single-cell partitioning technology of the BD Rhapsody™ System. This scATAC-Seq assay showed high sensitivity (many unique fragments of appropriate size per cell) and specificity (most fragments clustered into distinct peaks, with significant enrichment around annotated Transcription Start Sites).

Here we demonstrate the power of this multiomics assay for exploring the regulation of gene expression through chromatin accessibility. Cell types were annotated, and we found enrichment of transcription factor motifs within CARs in each distinct cell type. Gene accessibility correlated well with RNA expression within cell types, and we demonstrate that more in-depth examination of gene regions can indicate the presence of enhancer elements.

## Methods

### Workflow of Single-Cell ATAC-Seq and mRNA Whole Transcriptome Analysis on BD Rhapsody™ HT Xpress System



**Figure 1.** Following the isolation of nuclei from cells, they undergo a tagmentation process utilizing a custom Tn5 transposase. This critical step involves the selective cutting of accessible chromatin regions, during which adapter sequences (pre-loaded onto the Tn5) are concurrently attached to both ends of the resulting fragments. The tagmented nuclei are then precisely loaded into a BD Rhapsody™ 8-Lane Cartridge, where they are individually partitioned into wells equipped with BD Rhapsody™ Enhanced Cell Capture Beads. This process ensures the simultaneous capture of both ATAC and mRNA fragments on the beads. Subsequently, distinct Whole Transcriptome Amplification (WTA) and ATAC libraries are prepared from these beads. These libraries are then sequenced and processed through a comprehensive analysis pipeline, leading to the generation of high-quality data that is fully compatible with existing bioinformatics tools, ready for in-depth analysis.

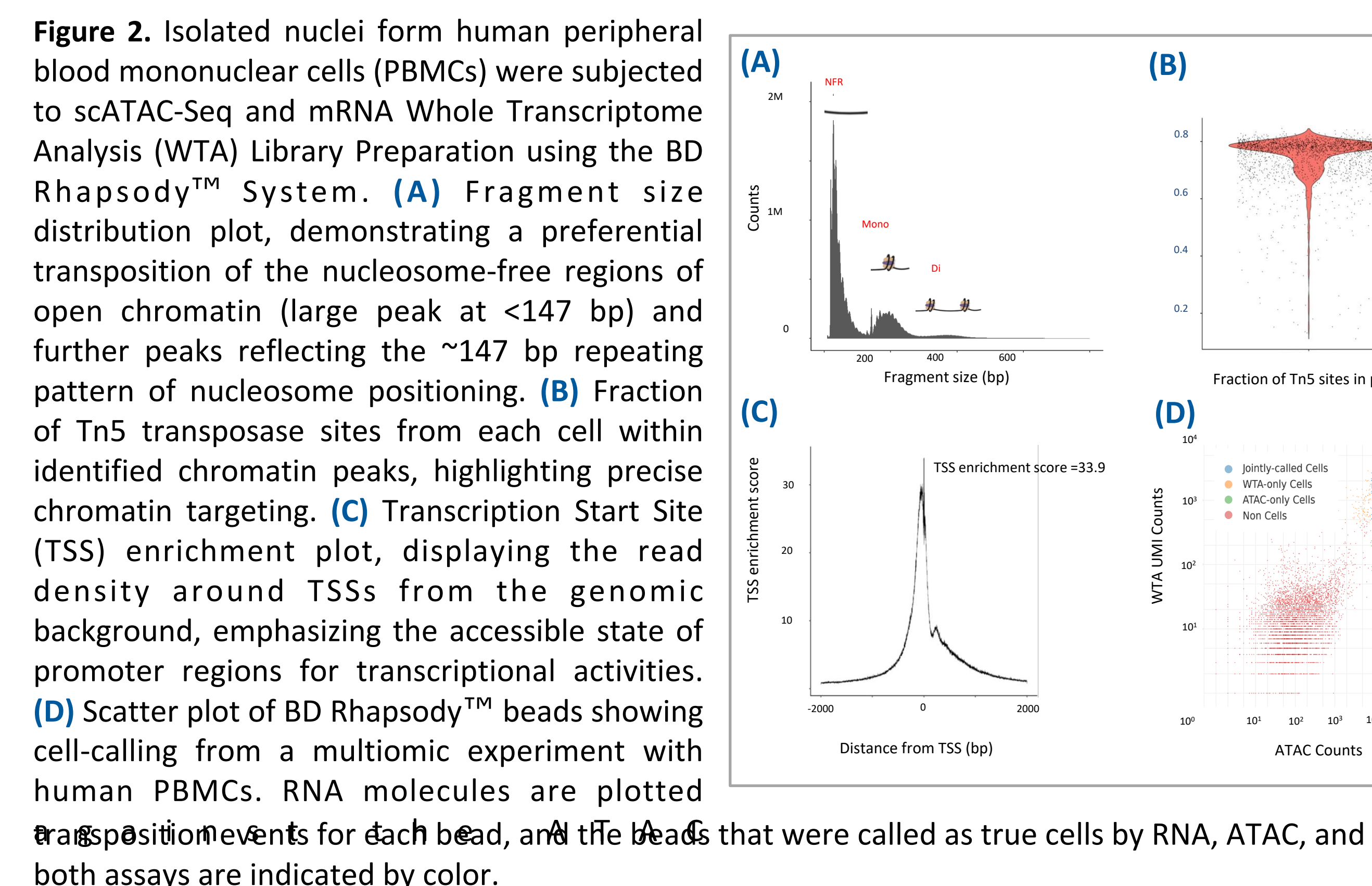
## Summary

**Empowering Molecular Biology and Immunology research with an accessible Multiomic scATAC-Seq Assay and analysis pipeline.**

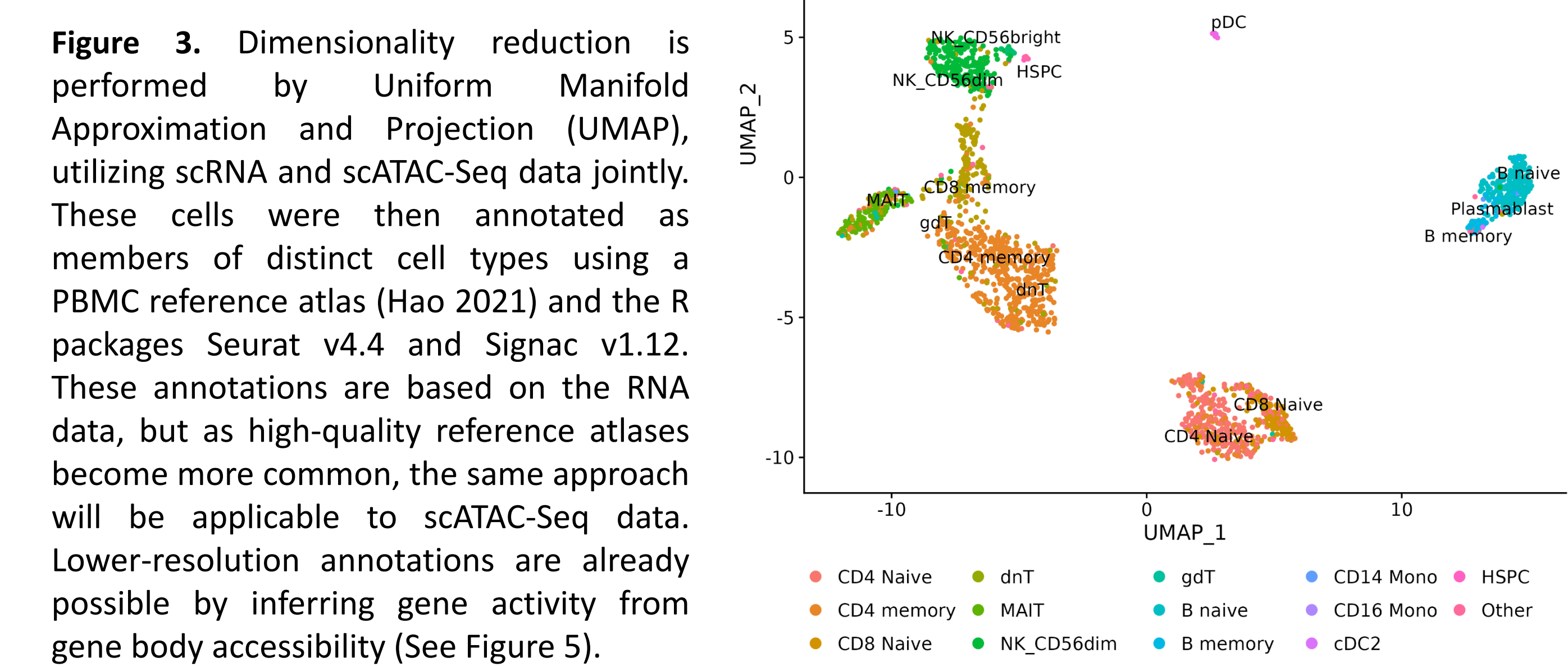
- Our study successfully integrates scRNA-Seq and scATAC-Seq through the BD Rhapsody™ Single-cell Analysis System for comprehensive multiomic analysis of PBMCs from multiple donors.
- This gentle approach allows for the simultaneous examination of gene expression and chromatin accessibility, providing a deeper understanding of cellular heterogeneity, development, and gene regulatory mechanisms.
- Our analysis reveals enriched transcription factor motifs and demonstrate a correlation between mRNA profiles and chromatin accessibility regions (CARs).
- Our data demonstrate consistency across both standalone scATAC-Seq and multiomic scATAC-Seq + scRNA-Seq analyses.
- Results of our pipeline are formatted for immediate compatibility with commonly used bioinformatic tools, allowing a quick and easy start to downstream analysis.

## Results

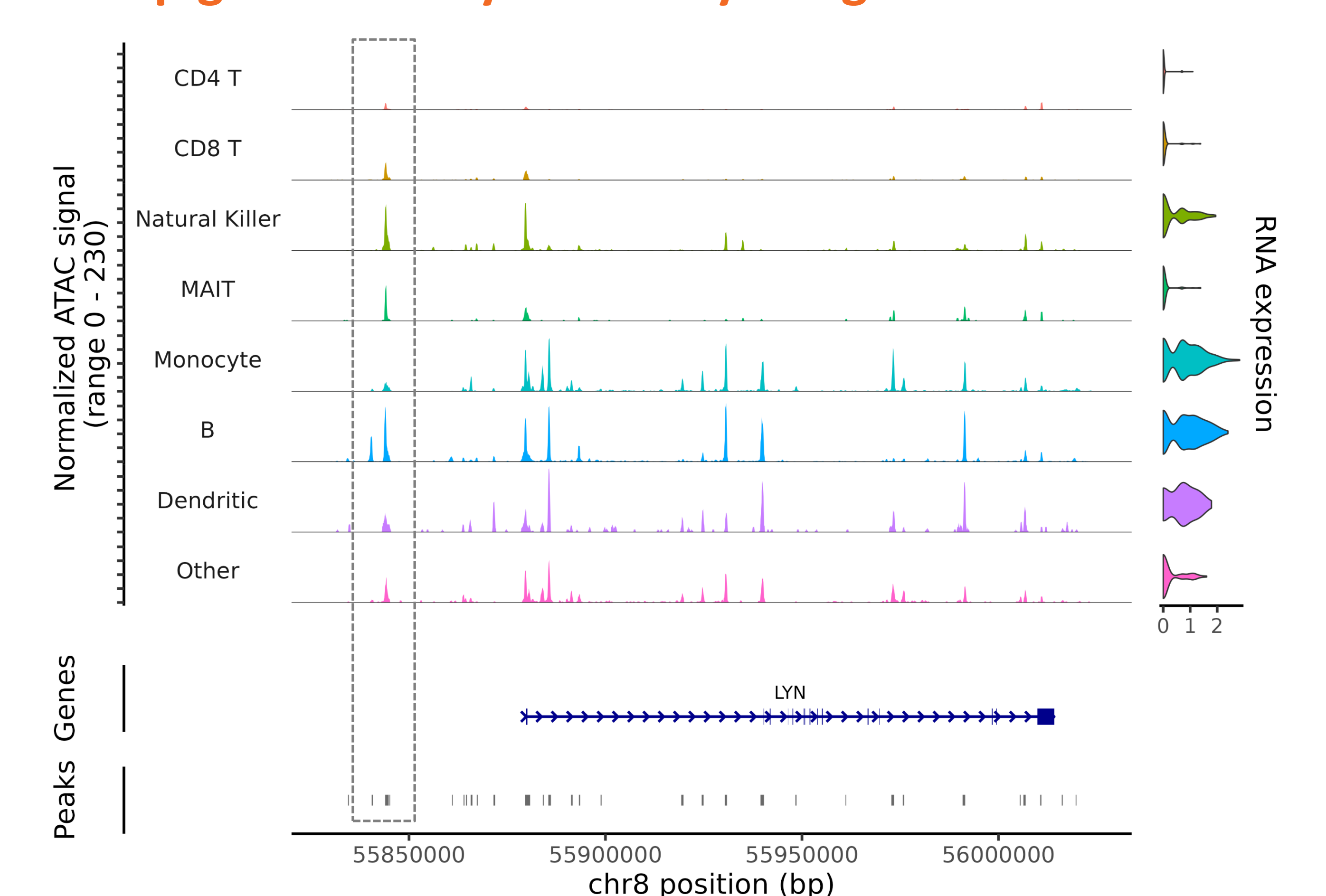
### Sensitivity and specificity metrics for the BD Rhapsody™ Single-Cell ATAC-Seq Assay



### RNA and ATAC data can be combined for dimensionality reduction

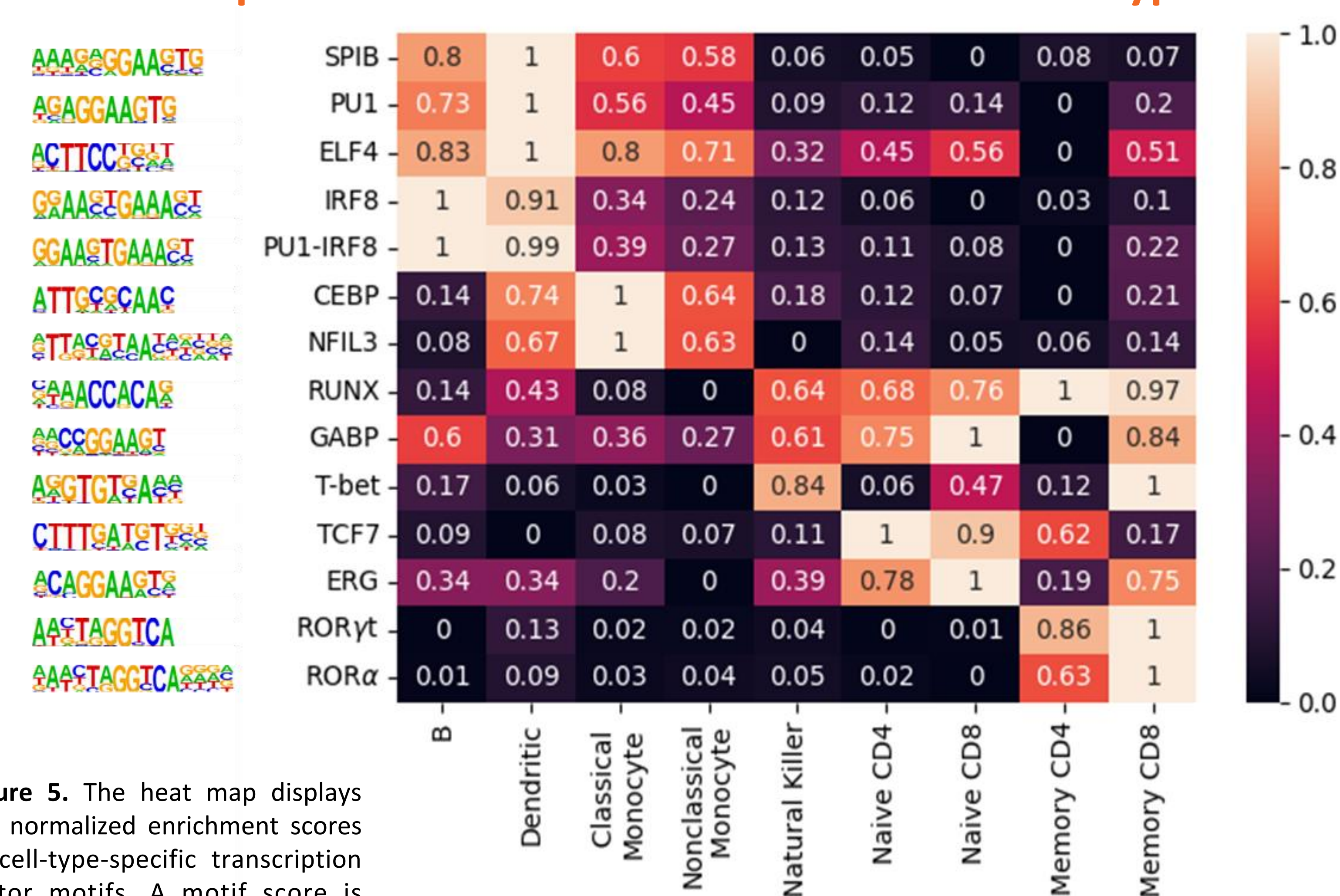


### Epigenome is dynamic beyond gene boundaries



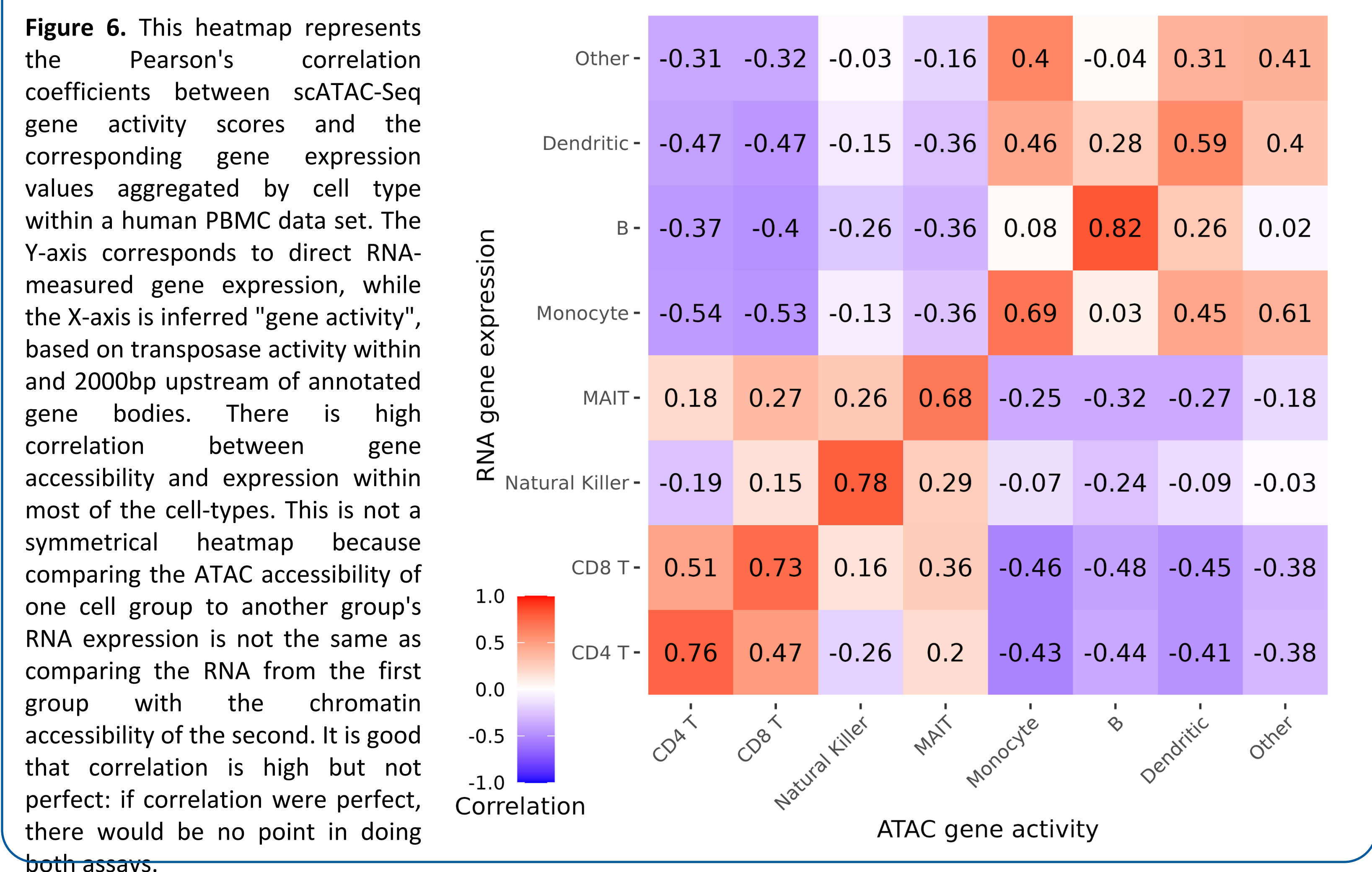
**Figure 4.** Aggregate ATAC signal within each cell type in the region surrounding the differentially expressed gene LYN. Log-normalized expression of LYN in the cells of each group is displayed in violin plots. Detected peaks of transposase activity are shown below the gene annotation. Note that some obvious peaks are visible upstream of the gene TSS, indicating a likely enhancer element (dotted box).

### Transcription factor motif enrichment across cell types



**Figure 5.** The heatmap displays the normalized enrichment scores of cell-type-specific transcription factor motifs. A motif score is calculated based on binomial distribution, using Homer for each pair of a cell type and a transcription factor, it calculates a relative motif enrichment in the differentially accessible regions of the given cell type, relative to randomly-selected background regions with controlling GC-content. Then the motif scores were normalized across cell types for each motif, with a scale ranging from 0 to 1. A score of 0 represents the least enrichment, while a score of 1 represents the highest enrichment of the motif. Models of transcription factor binding motifs are displayed to the left.

### Expression correlates with gene accessibility within cell types



### High reproducibility across PBMC replicates with BD Rhapsody™ multiomic scATAC-Seq assays: A peak comparison analysis

